

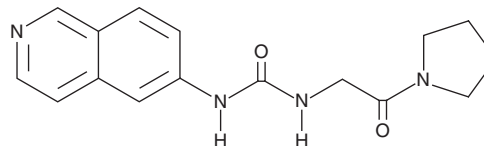
PRODUCT INFORMATION



SGC707

Item No. 17017

CAS Registry No.: 1687736-54-4
Formal Name: N-6-isoquinolinyl-N'-[2-oxo-2-(1-pyrrolidinyl)ethyl]-urea
MF: C₁₆H₁₈N₄O₂
FW: 298.3
Purity: ≥98%
UV/Vis.: λ_{max}: 248, 294 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SGC707 is supplied as a crystalline solid. A stock solution may be made by dissolving the SGC707 in the solvent of choice. SGC707 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SGC707 in ethanol is approximately 5 mg/ml and approximately 2 mg/ml in DMSO and DMF.

Description

Protein arginine N-methyltransferase 3 (PRMT3, Item No. 11642) is a predominantly cytoplasmic enzyme that is constitutively expressed.^{1,2} SGC707 is a potent allosteric inhibitor of PRMT3 (IC₅₀ = 50 nM) with >100-fold selectivity over other methyltransferases and other non-epigenetic targets. Developed by the Structural Genomics Consortium (SGC), SGC707 avidly binds PRMT3 (K_d = 50 nM by isothermal titration calorimetry) and inhibits the methylation of histones in cells with an IC₅₀ value below 1 μM.

References

1. Tang, J., Gary, J.D., Clarke, S., *et al.* PRMT3, a type I protein arginine N-methyltransferase that differs from PRMT1 in its oligomerization, subcellular localization, substrate specificity, and regulation. *J. Biol. Chem.* **273**(27), 16935-16945 (1998).
2. Wolf, S.S. The protein arginine methyltransferase family: An update about function, new perspectives and the physiological role in humans. *Cell. Mol. Life Sci.* **66**, 2109-2121 (2009).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/14/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM